

## Rose- $\beta$ -D-Gal

Rose- $\beta$ -D-Gal (6-Chloro-3-indoxyl- $\beta$ -D-galactopyranoside) is a chromogenic substrate for  $\beta$ -galactosidase. The product is similar to X-gal, but generates an intense pink colored precipitate ( $\lambda_{\max} \sim 540$  nm) on enzymatic hydrolysis.

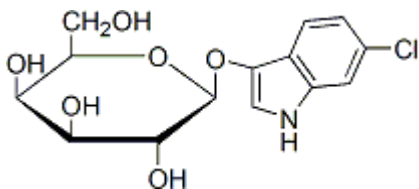


### Product attributes

## Product Description

Rose- $\beta$ -D-Gal (6-Chloro-3-indoxyl- $\beta$ -D-galactopyranoside) is a chromogenic substrate for  $\beta$ -galactosidase. The product is similar to X-gal, but generates an intense pink colored precipitate ( $\lambda_{\max} \sim 540$  nm) on enzymatic hydrolysis.

- White solid soluble in DMSO
- Store at 4 °C and protect from light
- MW: 329.74
- C<sub>14</sub>H<sub>16</sub>ClNO<sub>6</sub>



## References

1. Molecular Cloning: A laboratory Manual, Cold Spring Harbor Laboratory, B. 14, p186 (1989).
2. Biotechniques 27, 438 (1999).
3. Oncogene 10, 2323 (1995).
4. Dev Biol 161, 77 (1994).
5. J Histochem Cytochem 42, 1299 (1994).
6. BioTechniques 15, 974 (1993).
7. BioTechniques, 15, 292 (1993).

This datasheet was generated on June 4, 2026 at 08:01:05 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/rose-b-d-gal-6-chloro-3-indoxyl-b-d-galactopyranoside/>